

### Human TIMP1 ELISA Kit ab100651

[6 References](#) [2 图像](#)

#### 概述

产品名称	人TIMP1 ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 40 pg/ml
范围	24.69 pg/ml - 18000 pg/ml
回收率	93 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	94.49	84% - 104%
Serum	92.43	81% - 102%
Plasma	93.67	83% - 102%

**实验步骤** Multiple steps standard assay

**种属反应性** 与反应: Human

**产品概述** Abcam's TIMP1 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human TIMP1 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Human TIMP1 coated on a 96-well plate. Standards and samples are pipetted into the wells and TIMP1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human TIMP1 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of TIMP1 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Get higher sensitivity in only 90 minutes with Human TIMP1 ELISA Kit ([ab187394](#)) from our SimpleStep ELISA<sup>®</sup> range.

**说明** Optimization may be required with urine samples.

**平台** Microplate

## 性能

**存放说明** Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human TIMP1	2 vials
Recombinant Human TIMP1 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TIMP1 Microplate (12 x 8 wells)	1 unit
TMB One-Step Substrate Reagent	1 x 12ml

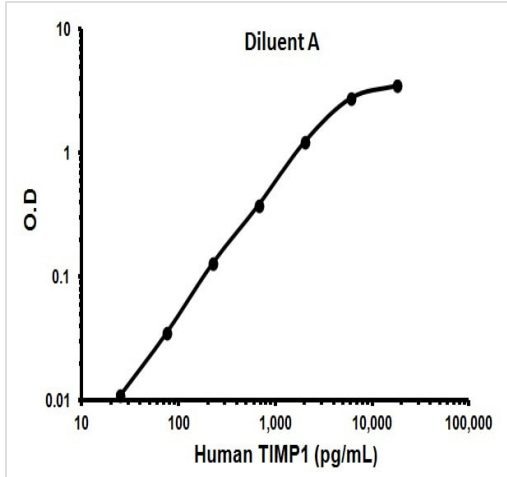
**功能** Complexes with metalloproteinases (such as collagenases) and irreversibly inactivates them by binding to their catalytic zinc cofactor. Also mediates erythropoiesis in vitro; but, unlike IL-3, it is species-specific, stimulating the growth and differentiation of only human and murine erythroid progenitors. Known to act on MMP-1, MMP-2, MMP-3, MMP-7, MMP-8, MMP-9, MMP-10, MMP-11, MMP-12, MMP-13 and MMP-16. Does not act on MMP-14.

**序列相似性** Belongs to the protease inhibitor I35 (TIMP) family.  
Contains 1 NTR domain.

**翻译后修饰** The activity of TIMP1 is dependent on the presence of disulfide bonds.

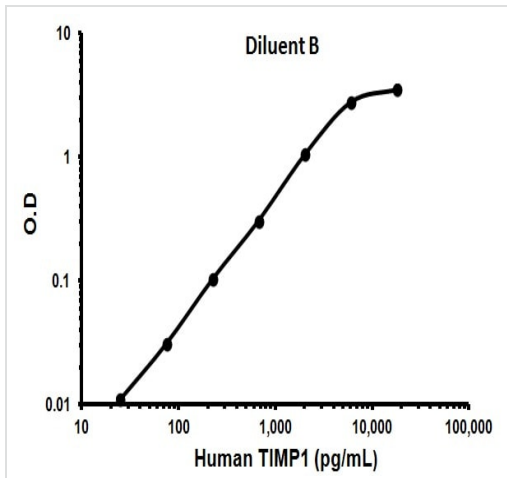
**细胞定位** Secreted.

## 图片



Representative standard curve using ab100651

Typical standard curve



Representative standard curve using ab100651

Typical standard curve

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

**Our Abpromise to you: Quality guaranteed and expert technical support**

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours

- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.cn/abpromise> or contact our technical team.

### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors